

## Rabbit Anti-Alexa Fluor 647 Antibody

Catalog Number: 46560

Unit Size: 1 mg

### Product Details

---

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
--------------------	--

Expiration Date	12 months upon receiving
-----------------	--------------------------

### Chemical Properties

---

Appearance	Liquid
------------	--------

Molecular Weight	N/A
------------------	-----

Soluble In	PBS
------------	-----

### Spectral Properties

---

Excitation Wavelength	N/A
-----------------------	-----

Emission Wavelength	N/A
---------------------	-----

### Applications

---

Rabbit Anti-Alexa Fluor 647 Antibody is a polyclonal antibody that specifically binds to Alexa Fluor® 647-conjugated molecules. It is ideal for detecting Alexa Fluor® 647-labeled antibodies, proteins, or probes.

The antibody may cross-react with related fluorophores such as iFluor® 647, Cy5, ATTO 647 and DyLight® 650. It is ideal for applications requiring signal amplification or dye conjugation verification.

*Alexa Fluor® and DyLight® are registered trademarks of Thermo Fisher Scientific.*